

D 2 Hydroxyglutarate Assay Kit (Colorimetric) ab211070

★★★★★ [1 Abreviews](#) [13 References](#) [2 Images](#)

Overview

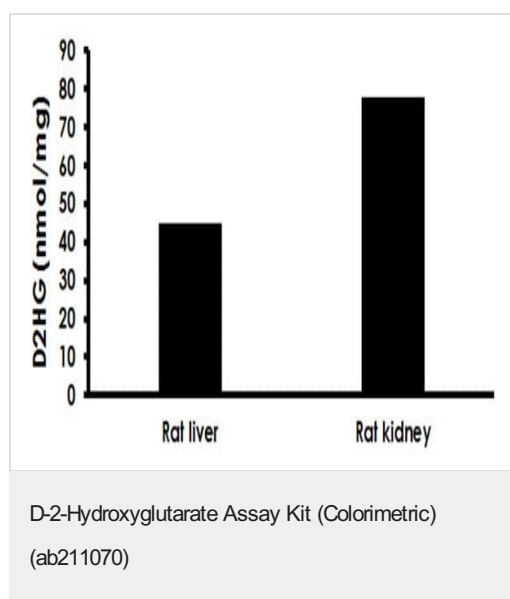
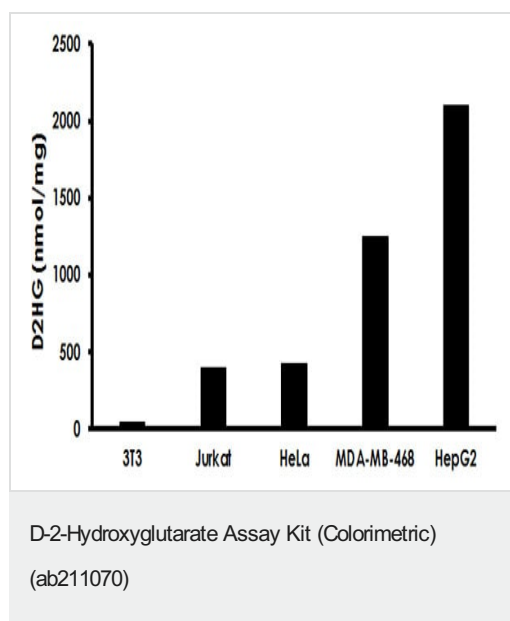
Product name	D 2 Hydroxyglutarate Assay Kit (Colorimetric)
Detection method	Colorimetric
Sample type	Urine, Serum, Plasma, Cell Lysate, Tissue Lysate
Assay type	Quantitative
Sensitivity	10 µM
Species reactivity	Reacts with: Mammals, Other species
Product overview	D-2-Hydroxyglutarate Assay Kit (ab211070) provides a convenient method to detect D-2-Hydroxyglutarate (D2HG) in urine, cell and tissue lysates. In this assay, D-2-Hydroxyglutarate is oxidized to α-Ketoglutarate in the presence of D2HG Enzyme and Substrate Mix. The intermediate reduces the probe to a colored product with strong absorbance at OD 450 nm. This absorbance is proportional to the amount of D2HG present in the samples. This assay kit is fast, sensitive, easy to use and high-through adaptable. It can measure D-2-Hydroxyglutarate levels less than 10 µM in various samples. This product is specific for D-2-Hydroxyglutarate and does not detect L-2-Hydroxyglutarate.
Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K213 D-2-Hydroxyglutarate (D2HG) Assay Kit (Colorimetric). K213-100 is the same size as the 100 test size of ab211070.</p> <p>In eukaryotic cells, Isocitrate Dehydrogenase (IDH1, IDH2 and IDH3) catalyzes the interconversion of Isocitrate and α-Ketoglutarate. In human cancers, an IDH mutation causes a gain-of-function, which reduces its affinity for isocitrate and facilitates the conversion of αketoglutarate to D-2-Hydroxyglutarate in the presence of NADP. D-2-Hydroxyglutarate (D2HG, D-2-Hydroxyglutarate) is present at low level in normal cells and tissues, but is significantly elevated in metabolic diseases and various cancers. Therefore, detection of elevated D2HG is important for early diagnosis, prognosis and the development of therapeutic strategies against these maladies.</p>
Platform	Microplate reader

Properties

Storage instructions	Store at -20°C. Please refer to protocols.
-----------------------------	--

Components	100 tests
Assay Buffer XXXII	1 x 20ml
D2HG Enzyme	1 vial
D2HG Standard	1 vial
Developer Solution III	1 vial

Images



Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours

- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit [**https://www.abcam.com/abpromise**](https://www.abcam.com/abpromise) or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors